

MA150C Moisture Analyzer

Fully Automatic Infrared Moisture Analyzer with Ceramic Heating Element

Moisture Analyzer



- Time-saving and convenient alternative to a drying oven
- Utmost measurement accuracy and durability "made in Germany"
- 20 program memory modules for quick method changing
- Very even heat distribution, ideally suited for thermally sensitive samples
- Optional version without glass, for use in food production applications

Max. weighing capacity	150 g
Repeatability, typical	Starting at an initial sample weight of approx. 1 g: $\pm 0.2\%$ starting at an initial sample weight of approx. 5 g: $\pm 0.05\%$
Readability	1 mg, 0.01%
Typical sample quantity	5 – 15 g
Reading display	% moisture content % dry weight RATIO value g residue g/kg residue g/l residue
Temperature range and settings	40°C–180°C, stand-by temperature selectable from 40–100°C in increments of 1 degree
Sample heating	Infrared heating using a round ceramic heating element
Heating programs	Standard drying, gentle drying
Shutoff criteria	Optional: – Fully automatic – Semi-automatic (1–50 mg/5–300 sec.) – Semi-automatic (0.1–5.0%/5–300 sec.) – Timer setting (0.1–99.9 min.) – Manual
Access to sample chamber	Hinged cover
Measuring program	20 programs saved in a non-volatile memory (freely selectable parameters)
Data storage	Results are saved until the start of the next measurement
Operator guidance features	Menu-driven, alphanumeric dialogue text (5 languages selectable) and symbols
Parameter data entry	Via softkeys, menu-specific key allocation
Report printout	– Printout using the optional, external printer YDP20-OCE – Configurable GLP-report for measuring results calibration adjustment (5 languages selectable)



Open sample chamber



MA150C featuring the accessory YDS04MA:
To replace glass panels with metal panels for use
in food production

Test equipment monitoring	<ul style="list-style-type: none"> – External calibration using optional weight sets – Automatic temperature calibration with optional temperature adjustment set YTM03MA
Safety	Parameter settings password-protected against unauthorized access
Voltage frequency	48-60 Hz
Power consumption	Max. 700 VA
Operating temperature range	10°C...30°C
Housing dimensions (W × D × H)	213 × 320 × 180.5
Weight	Approx. 5.5 kg

Available Models

	Supply voltage
MA150C-000115V1	115 V +15%, -20%
MA150C-000230V1	230 V +15%, -20%

Accessories

6965542	Disposable sample pans, 80 units, aluminum, round, Ø 90 mm
6906940	Fiberglass pad, Ø 90 mm for pasty and fatty samples, hard quality, 80 units
6906941	Fiberglass pad, Ø 90 mm for liquid and fatty samples, soft quality, with high suction force, 200 units
YAT01MA	Single-use pipettes for weighing liquid samples, 500 units
YDS04MA	Panel replacement set, aluminum panels for replacing glass panels, to meet FDA HACCP regulations (conversion kit)
YSC02	SartoCollect, software for data communication between moisture analyzer and PC (connecting cable must be ordered separately) Prerequisite: MS Windows 2000 Windows XP Professional, Vista Ultimate, Windows 7 Ultimate.
YCW5128-02	External calibration weight, 100 g (E2) with DKD certificate
YTM03MA	Temperature adjustment set with manufacturer certificate for adjustment of the heat source
YDP20-OCE	Data printer, suitable for external verification, with date, time, statistics, numerator including cable
6906918	Ink ribbon for data printer
6906937	Paper rolls for data printer, 5 rolls, 50 m each

Equipment Supplied

- Moisture analyzer with data interface
- Power cord
- User manual
- 80 alu sample pans
- In-use dust cover for keypad
- 1 pair of forceps

Sartorius Weighing Technology GmbH
Weender Landstrasse 94–108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

info.mechatronics@sartorius.com
www.sartorius-mechatronics.com

Subject to change without notice.
Printed in Germany on paper that has been bleached
without any use of chlorine. | W
Publication No.: W--2033-e11102
Order No.: 98649-012-38